

# c[A(2',5')pG(3',5')p] / 3'2'-cGAMP

c[A(2',5')pG(3',5')p] is a cyclic dinucleotide structurally related to the metazoan second messenger and STING activator c[G(2',5')pA(3',5')p] (2'3'-cGAMP, Cat. No. C 161).

*BLG-C238-05 (5 x 1 µmol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.*

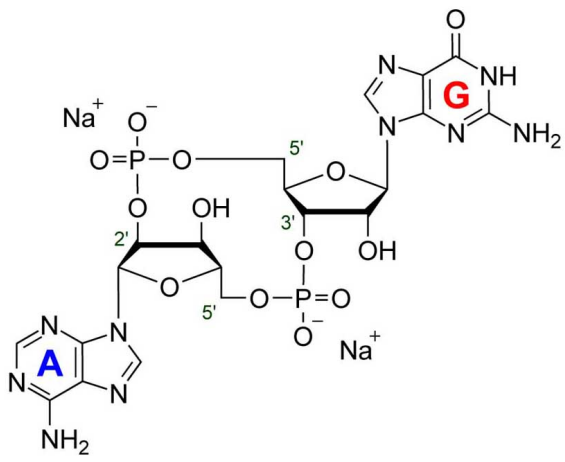
## Ordering Information

[Order Online »](#)

BLG-C238-01	1µmol
BLG-C238-05	5x1µmol

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Cyclic (adenosine-(2'→5')-monophosphate-guanosine-(3'→5')-monophosphate) . sodium salt, 3',5'-2',5'-cGAMP

**Couple Type** Modified nucleotides

**Formula**  $C_{20}H_{24}N_{10}O_{13}P_2$  (free acid)

**MW** 674.4 (free acid)

**Purity** > 98% HPLC

**Quantity** 0.5µmol is equivalent to ~337 µg

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)